

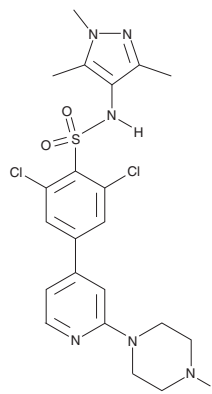
PRODUCT INFORMATION



DDD85646

Item No. 13839

CAS Registry No.: 1215010-55-1
Formal Name: 2,6-dichloro-4-[2-(1-piperazinyl)-4-pyridinyl]-N-(1,3,5-trimethyl-1H-pyrazol-4-yl)benzenesulfonamide
Synonym: IMP-366
MF: C₂₁H₂₄Cl₂N₆O₂S
FW: 495.4
Purity: ≥95%
UV/Vis.: λ_{max}: 226, 258 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

DDD85646 is supplied as a crystalline solid. A stock solution may be made by dissolving the DDD85646 in the solvent of choice. DDD85646 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of DDD85646 in these solvents is approximately 25 mg/ml. DDD85646 is also slightly soluble in ethanol.

DDD85646 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DDD85646 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. DDD85646 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

DDD85646 is a moderately bioavailable pyrazole sulphonamide inhibitor of *T. brucei* N-myristoyltransferase (TbNMT) with an apparent K_i value of 1.44 nM.¹ *T. brucei* is the parasite responsible for human African trypanosomiasis (HAT), also known as African sleeping sickness, and is transmitted through the tsetse fly. DDD85646 administration in mice (12.5 mg/kg for four days) ameliorated *T. brucei* in an acute mouse model of HAT.¹ It less potently inhibits growth of *T. cruzi* (EC₅₀ = 6.9 μM).

Reference

1. Frearson, J.A., Brand, S., McElroy, S.P., *et al.* N-myristoyltransferase inhibitors as new leads to treat sleeping sickness. *Nature* **464(7829)**, 728-732 (2010).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/04/2020

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM